

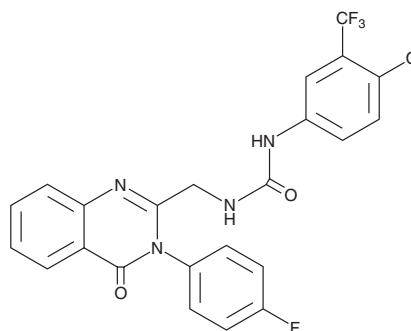
PRODUCT INFORMATION



Hedgehog Antagonist VIII

Item No. 16097

CAS Registry No.: 330796-24-2
Formal Name: N-[4-chloro-3-(trifluoromethyl)phenyl]-N'-[[3-(4-fluorophenyl)-3,4-dihydro-4-oxo-2-quinazolinyl]methyl]-urea
Synonym: Hh Antagonist VIII
MF: C₂₃H₁₅ClF₄N₄O₂
FW: 490.8
Purity: ≥98%
UV/Vis.: λ_{max}: 225, 253 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Hedgehog (Hh) antagonist VIII is supplied as a crystalline solid. A stock solution may be made by dissolving the Hh antagonist VIII in the solvent of choice, which should be purged with an inert gas. Hh antagonist VIII is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of Hh antagonist VIII in these solvents is approximately 10, 25, and 30 mg/ml, respectively.

Hh antagonist VIII is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, Hh antagonist VIII should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Hh antagonist VIII has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Hh proteins, important regulators of development, bind the cell-surface protein Patched, allowing activation of Smoothened. In vertebrates, this ultimately leads to the activation of the zinc-finger transcription factors of the Gli family. Overactivation of this pathway contributes to certain cancers, including glioblastoma, for which the Gli proteins are named. Hh antagonist VIII is a cell-permeable quinazolinyl-urea compound that has been shown to inhibit Gli transcription activity with an IC₅₀ value of 70 nM.¹

Reference

1. Brunton, S.A., Stibbard, J.H., Rubin, L.L., *et al.* Potent inhibitors of the hedgehog signaling pathway. *J. Med. Chem.* **51**(5), 1108-1110 (2008).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/11/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM